

# TREVIGEN<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

## Cultrex<sup>™</sup> Mouse Laminin I

Catalog #: 3400-010-01

Size: 1 mg

**Description:** This highly purified preparation of mouse laminin I increases cell adhesion, migration, growth, and differentiation. It is composed of  $\alpha 1\beta 1\gamma 1$  chains with a total Mr of 800,000 and is used for coating of culture dishes.

**Source:** Murine Engelbreth-Holm-Swarm (EHS) tumor.

**Storage Buffer:** Dulbecco's Modified Eagle's Medium with 10  $\mu$ g/ml gentamycin sulfate.

**Storage:** Store at -20°C or -80°C in a manual defrost freezer.

**Purity:** Purity >90% by SDS-PAGE.

### Specifications:

#### Functional Assays:

- Supports the formation of neuronal filaments of NG108-15 cells in a neurite outgrowth assay.

#### Sterility Testing:

- No bacterial or fungal growth detected after incubation at 37°C for 14 days following USP XXIV, Chapter 71 sterility testing.
- No mycoplasma contamination detected by PCR
- Endotoxin concentrations  $\leq$  20 EU/ml by LAL assay.

### Coating procedure:

The recommended working concentration is 0.05-10  $\mu$ g/cm<sup>2</sup> of growth surface (0.05 -10  $\mu$ g/ml) depending on cell type.

**A.** Thaw stock solution on ice for several hours. Place plates on ice and prechill pipette tips. In a laminar flow hood, dilute appropriately with cold serum-free, cell culture medium. Mix and transfer to the wells of cold tissue culture plates. Spread the solution to completely cover the bottom of the wells.

**B.** The following table is a guide for the suggested volumes required per well:

<u>Plate type</u>	<u>Volume Laminin per Well</u>
6 wells (or 35 mm dish)	1 ml
24 wells	200 $\mu$ l
48 wells	50 $\mu$ l
96 wells	20 $\mu$ l

**C.** Incubate the plates at 37°C for 1 hour. In the laminar flow hood, remove excess liquid from the wells of the tissue culture plate.

**D.** Rinse the wells once with tissue culture medium and then add your cells.

**Reference:** Kleinman H.K. 1999. The Laminins. *Proc. Indian Acad. Sci. (Chem.Sci.)*, 111:179-1841.



**Mouse Laminin I**

**Catalog #: 3400-010-01**

**Storage: -20°C or -80°C**

**TREVIGEN<sup>®</sup>**

**1-800-873-8443**

**TREVIGEN<sup>®</sup>**

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com